

Separation of some methylated phenols

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F254s (P/N 1. 13192, alt P/N 1. 15647)
Eluent	Chloroform/methanol 95/5 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. p-Cresol 2. 2,3-Dimethylphenol 3. 2,4,6-Trimethylphenol (all 0,2%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

